

## WASP (Ab-290) Antibody [#B0597]

<b>Catalog Number:</b>	B0597
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P42768
<b>Other Names:</b>	IMD2, WAS, Wiskott-Aldrich syndrome protein, Wiskott-Aldrich syndrome protein homolog
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human WASP around the phosphorylation site of tyrosine 290 (L-I-Y <sup>P</sup> -D-F).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	WASP (Ab-290) antibody detects endogenous levels of total WASP protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Hervé Falet, Blood, Aug 2002; 100: 2113 - 2122. E Remold-O'Donnell, J. Immunol., May 1997; 158: 4021. Ornella Parolini, Blood, Jul 1997; 90: 70 - 75. Anna Shcherbina, J. Immunol., Dec 1999; 163: 6314.